

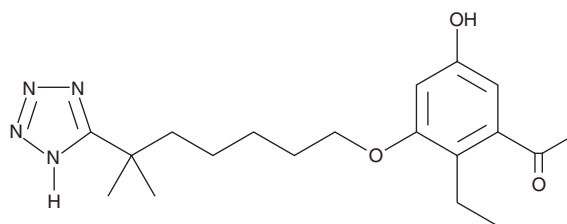
PRODUCT INFORMATION



LY255283

Item No. 70715

CAS Registry No.: 117690-79-6
Formal Name: 1-[5-ethyl-2-hydroxy-4-[[6-methyl-6-(1H-tetrazol-5-yl)heptyl]oxy]phenyl]-ethanone
MF: C₁₉H₂₈N₄O₃
FW: 360.5
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 232, 276, 326 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

LY255283 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY255283 in the solvent of choice, which should be purged with an inert gas. LY255283 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LY255283 in these solvents is approximately 1, 30, and 50 mg/ml, respectively.

LY255283 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY255283 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LY255283 has a solubility of approximately 100 µg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The tetrazole LY255283 is a competitive antagonist of the BLT₂ receptor. It displaces radiolabeled LTB₄ from guinea pig lung membrane, with an IC₅₀ of about 100 nM.¹ LY255283 exhibits IC₅₀ values of ~950 nM and >10 µM at human recombinant BLT₂ and BLT₁ receptors, respectively.² LY255283 inhibits eosinophil chemotaxis by 80% at a concentration of 10 µM, and inhibits the binding of radiolabeled LTB₄ to eosinophil membranes with an IC₅₀ of 260 nM.³

References

1. Silbaugh, S.A., Stengel, P.W., Cockerham, S.L., *et al.* Pulmonary actions of LY255283, a leukotriene B₄ receptor antagonist. *Eur. J. Pharmacol.* **223**, 57-64 (1992).
2. Yokomizo, T., Kato, K., Hagiya, H., *et al.* Hydroxyeicosanoids bind to and activate the low affinity leukotriene B₄ receptor, BLT₂. *J. Biol. Chem.* **276**(15), 12454-12459 (2001).
3. Richards, I.M., Sun, F.F., Taylor, B.M., *et al.* Contribution of leukotriene B₄ to airway inflammation and the effect of antagonists. *Ann. N. Y. Acad. Sci.* **629**, 274-287 (1991).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/05/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM